

beta Galactosidase Staining Kit ab102534

★★★★★ [1 Abreviews](#) [15 References](#) [画像数 1](#)

医薬用外劇物

製品の概要

製品名	beta Galactosidase Staining Kit
サンプルの種類	Adherent cells, Suspension cells
アッセイタイプ	Direct
製品の概要	<p>The LacZ gene from <i>E. coli</i> is one of the most commonly used reporter genes for testing the efficiency of expression vector mediated gene transfer and for studying the regulation of promoters of genes. The LacZ gene encodes the enzyme β-galactosidase, which is very stable, resistant to proteolytic degradation, can utilize a variety of substrates and can be easily assayed <i>in situ</i>.</p> <p>Abcam's beta Galactosidase Detection Kit (ab102534) utilizes X-gal as the substrate for β-galactosidase.</p> <p>Visit our FAQs page for tips and troubleshooting.</p>

特記事項

This product is manufactured by BioVision, an Abcam company and was previously called K802 Beta-Galactosidase Staining Kit. K802-250 is the same size as the 250 test size of ab102534.

This kit has been optimized for transfected cells. If you want to detect senescence, we recommend using [Senescence Detection Kit \(ab65351\)](#).

製品の特性

保存方法 Store at -20°C. Please refer to protocols.

内容	250 tests
Fixative Solution III	1 x 125ml
Staining Solution II	1 x 125ml
100X Staining Supplement	1 x 1.5ml
X-Gal	1 x 150mg

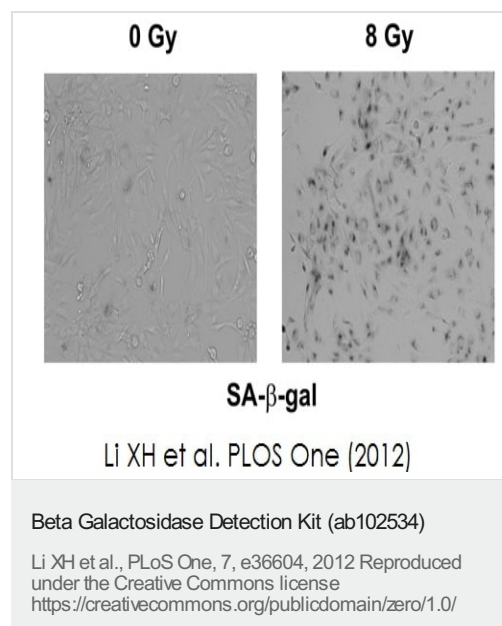
関連性 Beta galactosidase is a hydrolase enzyme that cleaves beta-linked terminal galactosyl residues

from gangliosides, glycoproteins, and glycosaminoglycans. Beta galactosidase is an essential enzyme in the human body. Deficiencies in the protein can result in galactosialidosis or Morquio B syndrome. Senescent cells display senescence-associated expression of beta galactosidase activity.

細胞内局在

Isoform 1: Lysosome. Isoform 2: Cytoplasm, perinuclear region. Note=Localized to the perinuclear area of the cytoplasm but not to lysosomes.

画像



Gamma-radiation induced senescence in hFOB cells. The number of senescence-associated-β-galactosidase positive cells (*left panel*) clearly increases 72h after 8 Gy γ irradiation. Representative data was from one of a total of 3 independent experiments. *Image obtained from Li XH et al; PLOS One; 2012;7(5):e36604*

Please note: All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.co.jp/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors